

Datasheet of GLuc-ON™ CREB Transcriptional Response Element (TRE) Lentiviral Clone (TR200)

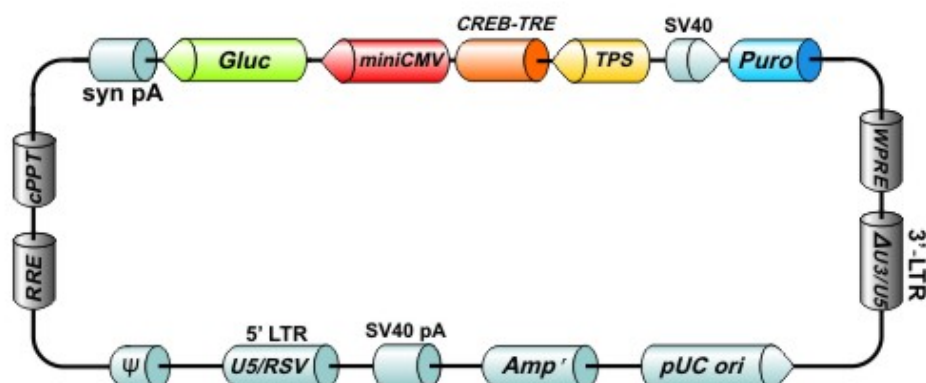


Figure 1. GLuc-ON™ CREB Transcriptional Response Element (TRE) Lentiviral Clone (TR200)

Catalog: TR200

Description:

The CREB transcriptional response element (TRE) lentiviral clone assays the status of the cAMP/PKA signaling pathway through the activity of the transcription factor CREB. The CREB transcriptional TRE clone contains a minimal promoter and tandem repeats of the CREB transcriptional response element upstream of a secreted Gaussia luciferase reporter gene. The number of response elements, as well as the intervening sequence between the elements has been optimized to maximize the sensitivity and signal to noise ratio.

Delivery: 500 ng/μL × 40 μL

Storage Conditions: -20°C

.Functional validation in common cell line

QC: .Full sequence verification

.Restriction Digestion

Validation Data

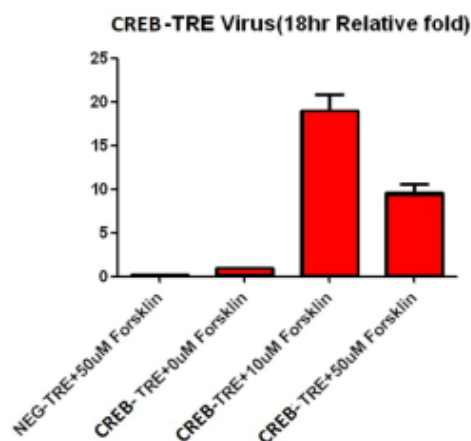


Figure 2. Validation data of GLuc-ON™ CREB Transcriptional Response Element (TRE) Lentiviral Clone (TR200).

Restriction Enzyme Information for TR200

Table 1. Restriction Enzymes That Do Not Cleave TR200

AfeI	AgeI	AscI	BclI	BlnI	BspEI	BsrGI	BstXI
BstZ17I	HpaI	MfeI	NdeI	NotI	PacI	PmeI	PshAI
SbfI	SgfI	SgrAI	SnaBI	SpeI	SrfI	Swal	XbaI

Table 2. Restriction Enzymes That Cleave TR200 Once

Acc16I	4220	AcvI	715	AjlI	942	AleI	6556	Alol	7848	Apal	516	AspI	1047	AsuII	8287
AviII	4220	AxyI	2339	BaeI	5338	BamHI	7879	BbrPI	715	BbvCI	6403	BmgBI	942	Bpu14I	8287
Bsa29I	8198	BsaAI	715	BsaBI	7840	Bse21I	2339	Bse8I	7840	BseCI	8198	BseJI	7840	BshVI	8198
BsiVI	1061	Bsp119I	8287	Bsp120I	512	BspDI	8198	BspT104I	8287	BstBAI	715	BstBI	8287	BstEII	1139
BstENI	6145	BstPI	1139	Bsu15I	8198	Bsu36I	2339	BtrI	942	Clal	8198	Cpol	1121	CspI	1121
Eco32I	8519	Eco72I	715	Eco81I	2339	Eco91I	1139	EcoNI	6145	EcoO65I	1139	EcoRI	8260	EcoRV	8519
FseI	6125	FspI	4220	KfiI	6913	MluI	1605	NbsI	4220	NspV	8287	OliI	6556	PaeR7I	7308
Pfi23II	1061	PfiFI	1047	PmaCI	715	PmlI	715	Ppu21I	715	PpuMI	6913	Psp5II	6913	PspCI	715
PspEI	1139	PspLI	1061	PspOMI	512	PspPPI	6913	PspXI	7308	PstI	2223	PsyI	1047	RigI	6125
Rsr2I	1121	RsrII	1121	Sfr274I	7308	SfuI	8287	SgrDI	4919	Slal	7308	Sspl	4802	StrI	7308
Tth111I	1047	XagI	6145	XcmI	5144	XhoI	7308								

Table 3. Restriction Enzymes That Cleave TR200 Twice

AanI	2684	AanI	4957	AarI	738	AarI	8124	AasI	3216	AasI	8055	Adel	762	Adel	1576
AflII	2532	AflII	5492	AhdI	3998	AhdI	7569	Ajul	7228	Ajul	8290	AspA2I	324	AspA2I	550
AspEI	3998	AspEI	7569	AssI	4478	AssI	8193	AsuNHI	2720	AsuNHI	8067	AvrII	324	AvrII	550
Ball	1380	Ball	7777	BauI	3281	BauI	4662	BcgI	1446	BcgI	4503	BfrI	2532	BfrI	5492
BinI	324	BinI	550	BmcAI	4478	BmcAI	8193	BmeRI	3998	BmeRI	7569	BmtI	2724	BmtI	8071
Bpu10I	6403	Bpu10I	8076	BsePI	1460	BsePI	5686	BsmBI	986	BsmBI	1397	Bsp68I	5808	Bsp68I	7765
BspHI	3825	BspHI	4833	BspOI	2724	BspOI	8071	BspQI	2992	BspQI	6077	BspTI	2532	BspTI	5492
BssHII	1460	BssHII	5686	BssSI	3281	BssSI	4662	Bst2BI	3281	Bst2BI	4662	BstAFI	2532	BstAFI	5492
BtuMI	5808	BtuMI	7765	CciI	3825	CciI	4833	Cfr9I	345	Cfr9I	1054	CsiI	91	CsiI	1572
DrallI	762	DrallI	1576	DrdI	3216	DrdI	8055	Dril	3998	Dril	7569	DseDI	3216	DseDI	8055
Eam1105I	3998	Eam1105I	7569	Eco147I	323	Eco147I	1336	EcoT22I	74	EcoT22I	146	Esp3I	986	Esp3I	1397
KroI	2130	KroI	6121	LguI	2992	LguI	6077	MabI	91	MabI	1572	MlsI	1380	MlsI	7777
MluNI	1380	MluNI	7777	Mph1103I	74	Mph1103I	146	MroNI	2130	MroNI	6121	MscI	1380	MscI	7777
Msp20I	1380	Msp20I	7777	MspCI	2532	MspCI	5492	NaeI	2132	NaeI	6123	NgoMIV	2130	NgoMIV	6121
NheI	2720	NheI	8067	NruI	5808	NruI	7765	NsiI	74	NsiI	146	PagI	3825	PagI	4833
Paul	1460	Paul	5686	PceI	323	PceI	1336	PciI	889	PciI	3108	PciSI	2992	PciSI	6077
PdiI	2132	PdiI	6123	PfoI	1480	PfoI	1891	PscI	889	PscI	3108	Psil	2684	Psil	4957
PteI	1460	PteI	5686	RcaI	3825	RcaI	4833	RruI	5808	RruI	7765	SapI	2992	SapI	6077
Scal	4478	Scal	8193	SexAI	91	SexAI	1572	SfiI	7968	SfiI	8093	Smal	347	Smal	1056
SseBI	323	SseBI	1336	StuI	323	StuI	1336	TspMI	345	TspMI	1054	Vha464I	2532	Vha464I	5492
XmaI	345	XmaI	1054	XmaJI	324	XmaJI	550	Zrml	4478	Zrml	8193	Zsp2I	74	Zsp2I	146

Table 4. Restriction Enzymes That Cleave TR200 Three Times

Acc36I	738	Acc36I	7394	Acc36I	8124	Acc65I	840	Acc65I	2345	Acc65I	8080	AcII	422	AcII	4224
AcII	4597	Asp718I	840	Asp718I	2345	Asp718I	8080	BciVI	1687	BciVI	3311	BciVI	4835	BfuAI	738
BfuAI	7394	BfuAI	8124	BfuI	1687	BfuI	3311	BfuI	4835	BpvUI	4368	BpvUI	5320	BpvUI	7538
BseX3I	1453	BseX3I	6119	BseX3I	6123	BspMI	738	BspMI	7394	BspMI	8124	BstZI	1453	BstZI	6119
BstZI	6123	Bsul	1687	Bsul	3311	Bsul	4835	Bvel	738	Bvel	7394	Bvel	8124	Cfr42I	1219
Cfr42I	2121	Cfr42I	7741	DinI	1214	DinI	1345	DinI	5614	EagI	1453	EagI	6119	EagI	6123
EclXI	1453	EclXI	6119	EclXI	6123	Eco52I	1453	Eco52I	6119	Eco52I	6123	Egel	1214	Egel	1345
Egel	5614	Ehel	1214	Ehel	1345	Ehel	5614	KasI	1212	KasI	1343	KasI	5612	Kpnl	844
Kpnl	2349	Kpnl	8084	Kspl	1219	Kspl	2121	Kspl	7741	Mly113I	1213	Mly113I	1344	Mly113I	5613
Mvrl	4368	Mvrl	5320	Mvrl	7538	NarI	1213	NarI	1344	NarI	5613	Ple19I	4368	Ple19I	5320
Ple19I	7538	PluTI	1216	PluTI	1347	PluTI	5616	Psp1406I	422	Psp1406I	4224	Psp1406I	4597	Pvul	4368
Pvul	5320	Pvul	7538	Pvull	2376	Pvull	2932	Pvull	8046	SacII	1219	SacII	2121	SacII	7741
Sfol	1214	Sfol	1345	Sfol	5614	Sfr303I	1219	Sfr303I	2121	Sfr303I	7741	SgrBI	1219	SgrBI	2121
SgrBI	7741	SspDI	1212	SspDI	1343	SspDI	5612								